



PRODUCT INFORMATION & MANUAL

Aspartate Aminotransferase Activity Assay Kit (Colorimetric) *NBP3-24469*

For research use only.
Not for diagnostic or therapeutic
procedures.

www.novusbio.com - P: 303.730.1950 - P: 888.506.6887 - F: 303.730.1966 - technical@novusbio.com

Novus kits are guaranteed for 6 months from date of receipt

Aspartate Aminotransferase Activity Assay Kit (Colorimetric)

Catalog No: NBP3-24469

Method: Colorimetric method

Specification: 96T (Can detect 43 samples without duplication)

Instrument: Microplate reader

Sensitivity: 1.1 IU/L

Detection range: 1.1-72.3 IU/L

Average intra-assay CV (%): 5.3

Average inter-assay CV (%): 6.8

- ▲ This kit is for research use only.
- ▲ Instructions should be followed strictly, changes of operation may result in unreliable results.
- ▲ Please kindly provide us the lot number (on the outside of the box) of the kit for more efficient service.

General information

▲ Intended use

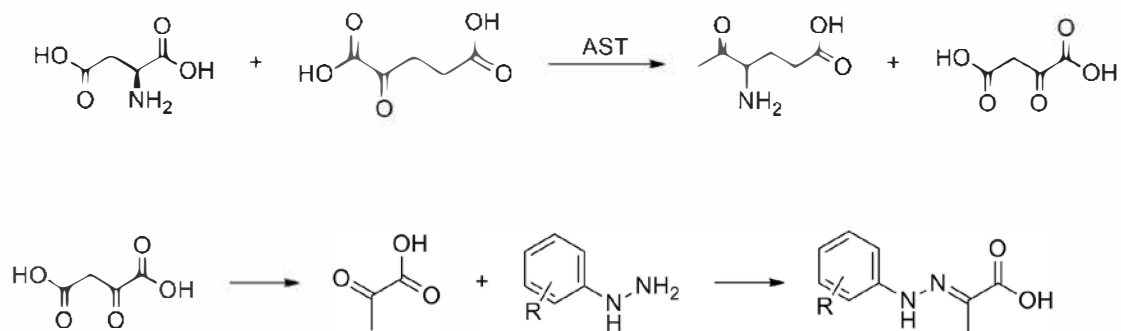
This kit can measure Aspartate Aminotransferase (AST/GOT) activity in animal serum (plasma), tissue, culture cells and cell culture supernatant, etc.

▲ Background

AST/GOT is a key enzyme in nitrogen metabolism, which is widely found in plasma and body tissues, including liver, heart, skeletal muscle, kidney, brain, pancreas, lung and erythrocyte. Changes in AST/GOT activity were found in acute pancreatitis, ischemic stroke, severe burns, periodontitis, acute renal disease and motor neuron disease.

▲ Detection principle

AST/GOT enables alpha-ketoglutaric acid and aspartic acid to displace amino and keto groups to form glutamic acid and oxaloacetic acid. Oxaloacetic acid can decarboxylate itself to form Pyruvic acid during the reaction. Pyruvic acid reacted with 2,4-dinitrophenylhydrazine (DNPH) to form 2,4-dinitrophenylhydrazone showing reddish brown in alkaline solution. Measure the OD values and calculate the enzyme activity.



▲ Kit components & storage

Item	Component	Specification	Storage
Reagent 1	Buffer Solution	0.5 mL × 1 vial	2-8°C , 12 months
Reagent 2	2 mmol/L Sodium Pyruvate	0.5 mL × 1 vial	2-8°C , 12 months
Reagent 3	Substrate Solution	5 mL × 1 vial	2-8°C , 12 months
Reagent 4	Chromogenic Agent	5 mL × 1 vial	2-8°C , 12 months, shading light
Reagent 5	Alkali Reagent	5 mL × 1 vial	2-8°C , 12 months
	Microplate	96 wells	No requirement
	Plate Sealer	2 pieces	

Note: The reagents must be stored strictly according to the preservation conditions in the above table. The reagents in different kits cannot be mixed with each other.

▲ Materials prepared by users

Instruments

Microplate reader (500-520 nm), Micropipettor, Vortex mixer, Incubator, Multichannel pipette

Reagents

Double distilled water, Normal saline (0.9% NaCl), PBS (0.01 M, pH 7.4)

▲ Safety data

Some of the reagents in the kit contain dangerous substances. It should be avoided to touch the skin and clothing. Wash immediately with plenty of water if touching it carelessly. All the samples and waste material should be treated according to the relevant rules of laboratory's biosafety.

▲ Precautions

Before the experiment, please read the instructions carefully, and wear gloves and work clothes.

▲ The key points of the assay

1. It is recommended to use multi-channel pipette to add reagent 5 working solution to reduce the difference between wells.
2. Detect the sample as soon as possible after collection. The serum sample can be store at 2-8°C for 7 days and -20°C for 20 days.

Pre-assay preparation

▲ Reagent preparation

1. Preparation of Reagent 5 working solution:

Dilute the Reagent 5 with double distilled water at the ratio of 1: 9 and mix fully. Prepare the fresh solution before use.

2. Incubate reagent 3 at 37 °C for 10 min.

▲ Sample preparation

The samples should be prepared as conventional methods. Also please refer to appendix II.

▲ Dilution of sample

It is recommended to take 2~3 samples with expected large difference to do pre-experiment before formal experiment and dilute the sample according to the result of the pre-experiment and the detection range (1.1-72.3 IU/L).

The recommended dilution factor for different samples is as follows (for reference only)

Sample type	Dilution factor
Human serum	1
Human plasma	1
Porcine serum	1
Rat serum	1
HC-60 cellular supernatant	1
Calu-3 cellular supernatant	1
10% Rat liver tissue homogenization	15-30
10% Rat lung tissue homogenization	2-8

Note: The diluent is normal saline (0.9% NaCl) or PBS (0.01 M, pH 7.4).

Assay protocol

▲ Plate set up

	1	2	3	4	5	6	7	8	9	10	11	12
A	A	A	S4	S4'	S12	S12'	S20	S20'	S28	S28'	S36	S36'
B	B	B	S5	S5'	S13	S13'	S21	S21'	S29	S29'	S37	S37'
C	C	C	S6	S6'	S14	S14'	S22	S22'	S30	S30'	S38	S38'
D	D	D	S7	S7'	S15	S15'	S23	S23'	S31	S31'	S39	S39'
E	E	E	S8	S8'	S16	S16'	S24	S24'	S32	S32'	S40	S40'
F	S1	S1'	S9	S9'	S17	S17'	S25	S25'	S33	S33'	S41	S41'
G	S2	S2'	S10	S10'	S18	S18'	S26	S26'	S34	S34'	S42	S42'
H	S3	S3'	S11	S11'	S19	S19'	S27	S27'	S35	S35'	S43	S43'

Note: A-E, standard wells; S1-S43, sample wells; S1'-S43, control wells.

▲ Detailed operating steps

(1) Standard wells: Add 5 μL of reagent 1 to the standard wells respectively (the pipettes should touch the bottom of the plate). Add 20, 18, 16, 14, 12 μL of reagent 3 to the standard wells from A to E, respectively. Add 0, 2, 4, 6, 8 μL of reagent 2 to the standard wells from A to E, respectively.

Sample wells: Add 20 μL of reagent 3 (pre-heated at 37°C for 10 min) and 5 μL of sample.

Control wells: Add 20 μL of reagent 3 (pre-heated at 37°C for 10 min).

(2) Mix fully (this is very important), then incubate at 37°C for 30 min.

(3) Add 20 μL of reagent 4 to each well.

- (4) Control wells: Add 5 μL of sample to Control wells.
- (5) Mix fully with microplate reader for 10 s, incubate at 37°C for 20 min.
- (6) Add 200 μL of reagent 5 working solution to each well (the multi-channel pipette is recommended).
- (7) Mix fully with microplate reader for 10 s, stand for 15 min at room temperature and measure the OD value of each well with microplate reader at 510 nm.

▲ Summary operation table

The preparation of standard curve

Set 5 wells of micro-plate for standard and operate according to the following operating table.

	A	B	C	D	E
Reagent 1 (μL)	5	5	5	5	5
Reagent 3 (μL)	20	18	16	14	12
Reagent 2 (μL)	0	2	4	6	8
Mix fully, then incubate at 37°C for 30 min.					
Reagent 4 (μL)	20	20	20	20	20
Mix fully and incubate at 37°C for 20 min.					
Reagent 5 working solution (μL)	200	200	200	200	200
Mix fully, stand for 15 min at room temperature and measure the OD value at 510 nm.					

The measurement of samples

	Control well	Sample well
Reagent 3 (μL) (pre-heated at 37°C for 10 min)	20	20
Sample (μL)		5
Mix fully, then incubate at 37°C for 30 min.		
Reagent 4 (μL)	20	20
Sample (μL)	5	
Mix fully and incubate at 37°C for 20 min.		
Reagent 5 working solution (μL)	200	200
Mix fully, stand for 15 min at room temperature and measure the OD value at 510 nm.		

▲ Calculation

1. Definition of international unit: The enzyme amount of $1\ \mu\text{mol}$ of NADH consumed in reaction system ($1\ \text{mL}$ sample or $1\ \text{g}$ tissue protein, 25°C) per minute is defined as 1 unit (wavelength is $340\ \text{nm}$, optical path is $1\ \text{cm}$).
2. Definition of Carmen unit: $1\ \text{mL}$ of sample, the total volume of reaction is $3\ \text{mL}$, wavelength is $340\ \text{nm}$, optical path is $1\ \text{cm}$, react at 25°C for $1\ \text{min}$, the amount of generated pyruvic acid which oxidize NADH to NAD^+ and cause absorbance decreasing 0.001 is as 1 unit. ($1\ \text{Carmen unit} = 0.482\ \text{IU/L}$, 25°C).

AST/GOT

3. Plot the standard curve by using OD value of standard and correspondent Carmen unit (0, 24, 61, 114, 190 Carmen unit) as x-axis and y-axis respectively. Create the standard curve with graph software (or EXCEL). The Carmen unit of the sample can be calculated according to the formula based on the OD value of sample.

The standard curve is $y=ax^2+bx+c$.

(1) Serum/plasma

$$\text{AST/GOT activity (IU/L)}=[a\times(\Delta A_{510})^2+b\times\Delta A_{510}+c]\times 0.482\times f$$

(2) Tissue and Cells

$$\text{AST/GOT activity (IU/gprot)}=[a\times(\Delta A_{510})^2+b\times\Delta A_{510}+c]\times 0.482\times f\div \text{Cpr}$$

Note:

y: Carmen unit.

x: $\text{OD}_{\text{standard}}-\text{OD}_{\text{blank}}$ (OD_{blank} is the OD value when the carmen unit is 0)

ΔA_{510} : $\text{OD}_{\text{sample}}-\text{OD}_{\text{control}}$

a, b, c: the constant of standard curve.

f: dilution factor of sample before tested.

Cpr: concentration of protein in sample (gprot/L)

Appendix I Data

▲ Example analysis

Take 5 μL of human serum, carry the assay according to the operation table.

The results are as follows:

standard curve: $y=2517.55x^2+74.50x+1.8995$, the average OD value of the sample is 0.259, the average OD value of the control is 0.233, and the calculation result is:

$$\text{AST activity (IU/L)} = [2517.55 \times (0.259 - 0.233)^2 + 74.5 \times (0.259 - 0.233) + 1.8995] \times 0.482 = 2.67 \text{ IU/L}$$

Appendix II Sample preparation

The following sample pretreatment methods are for reference only.

▲ Serum

Collect fresh blood and stand at 25°C for 30 min to clot the blood. Then centrifuge at 2000 g for 15 min at 4°C. Take the serum (which is the upper light yellow clarified liquid layer) to preserve it on ice for detection.

▲ Plasma

Take fresh blood into the tube which has anticoagulant (heparin is recommended), centrifuge at 700-1000 g for 10 min at 4°C. Take the plasma (which is the upper light yellow clarified liquid layer, don't take white blood cells and platelets in the middle layer) to preserve it on ice for detection.

▲ Urine

Collect fresh urine and centrifuge at 10000 g for 15 min at 4°C. Take the supernatant to preserve it on ice for detection.

▲ Tissue sample

Take 0.02-1g fresh tissue to wash with homogenization medium at 2-8°C. Absorb the water with filter paper and weigh. Homogenize at the ratio of the volume of homogenized medium (2-8°C) (mL): the weight of the tissue (g) =9:1, then centrifuge the tissue homogenate for 10 min at 1500 g at 4°C. Take the supernatant to preserve it on ice for detection. Meanwhile, determine the protein concentration of supernatant.

▲ Cells

Collect the cells and wash the cells with homogenization medium for 1~2 times. Centrifuge at 1000 g for 10 min and then discard the supernatant and keep the cell sediment. Add homogenization medium at a ratio of cell number (10^6): homogenization medium (μL) = 1: 300-500. Sonicate or grind with hand-operated in ice water bath. Centrifuge at 10000 g for 10 min, then take the supernatant and preserve it on ice for detection. Meanwhile, determine the protein concentration of supernatant.

Note:

1. Homogenized medium: PBS (0.01 M, pH 7.4) including 0.1 mM EDTA.
2. Homogenized method:
 - (1) Hand-operated: Weigh the tissue and mince to small pieces (1 mm^3), then put the tissues pieces to glass homogenized tube. Add homogenized medium into homogenized tube, place the tube into the ice bath with left hand, and insert the glass tamping rod vertically into the homogenized tube with the right hand to grind up and down for 6-8 min.
Or put the tissue into the mortar, and add liquid nitrogen to grind fully. Then add the homogenized medium to homogenize.
 - (2) Mechanical homogenate: Weigh the tissue to EP tube, add the homogenized medium to homogenize the tissue with homogenizer instrument (60 Hz, 90s) in the ice bath. (For samples of skin, muscle and plant tissue, the time of homogenization can be properly prolonged.)
 - (3) Ultrasonication: Treat the cells with ultrasonic cell disruptor (200 W, 2 s/ time, interval for 3 s, the total time is 5 min).

Appendix III References

1. Karmen A, Wroblewski F, Ladue J S. Transaminase activity in human blood. *Journal of Clinical Investigation*, 1995, 34(1): 126-131.
2. Desai S N, Desai P V. Aspartate aminotransferase and alanine aminotransferase activities of rat brain during crush syndrome. *Neuroscience Letters*, 2008, 447(1): 58-61.
3. Campos F, Sobrino T, Ramoscabrer P, et al. Neuroprotection by glutamate oxaloacetate transaminase in ischemic stroke: an experimental study. *J Cereb Blood Flow Metab*, 2011, 31(6): 1378-1386.
4. Hugon J, Tabaraud F, Rigaud M, et al. Glutamate dehydrogenase and aspartate aminotransferase in leukocytes of patients with motor neuron disease. *Neurology*, 1989, 39(7): 956-958.